

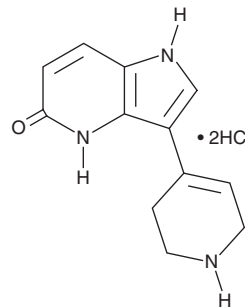
PRODUCT INFORMATION



CP 93,129 (hydrochloride)

Item No. 34548

CAS Registry No.: 879089-64-2
Formal Name: 1,4-dihydro-3-(1,2,3,6-tetrahydro-4-pyridinyl)-5H-pyrrolo[3,2-b]pyridin-5-one, dihydrochloride
MF: C₁₂H₁₃N₃O • 2HCl
FW: 288.2
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

CP 93,129 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the CP 93,129 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. CP 93,129 (hydrochloride) is soluble in the organic solvent DMSO. It is also soluble in water. The solubility of CP 93,129 (hydrochloride) in DMSO and water is approximately 100 mM. We do not recommend storing the aqueous solution for more than one day.

Description

CP 93,129 is an agonist of the serotonin (5-HT) receptor subtype 5-HT_{1B}.¹ It selectively binds 5-HT_{1B} over 5-HT_{1A}, 5-HT_{1C}, 5-HT_{1D}, and 5-HT₂ receptors (K_is = 0.0081, 1.5, 2.9, 1.1, and 7.2 μM, respectively). CP 93,129 (5.4 μM) inhibits potassium-induced GABA release from rat globus pallidus slices.² It reverses akinesia in a rat model of Parkinson's disease induced by reserpine (Item No. 16474) when administered intrapallidally at doses of 110 and 220 nmol/animal. CP 93,129 reduces food intake in rats when infused into the pons parabrachial nucleus or the paraventricular nucleus (PVN) of the hypothalamus (ED₅₀s = 1 and 50 nmol/animal, respectively).³ It also reduces the frequency of pursuits, tail rattles, threats, and attack bites in a resident-intruder test in mice.⁴

References

1. Koe, B.K., Nielsen, J.A., Macor, J.E., *et al.* Biochemical and behavioral studies of the 5-HT_{1B} receptor agonist, CP-94,253. *Drug Dev. Res.* **26(3)**, 241-250 (1992).
2. Chadha, A., Sur, C., Atack, J., *et al.* The 5HT_{1B} receptor agonist, CP-93129, inhibits [³H]-GABA release from rat globus pallidus slices and reverses akinesia following intrapallidal injection in the reserpine-treated rat. *Br. J. Pharmacol.* **130(8)**, 1927-1932 (2000).
3. Lee, M.D., Aloyo, V.J., Fluharty, S.J., *et al.* Infusion of the serotonin_{1B} (5-HT_{1B}) agonist CP-93,129 into the parabrachial nucleus potently and selectively reduces food intake in rats. *Psychopharmacology (Berl.)* **136(3)**, 304-307 (1997).
4. Bannai, M., Fish, E.W., Faccidomo, S., *et al.* Anti-aggressive effects of agonists at 5-HT_{1B} receptors in the dorsal raphe nucleus of mice. *Psychopharmacology (Berl.)* **193(2)**, 295-304 (2007).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/07/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM